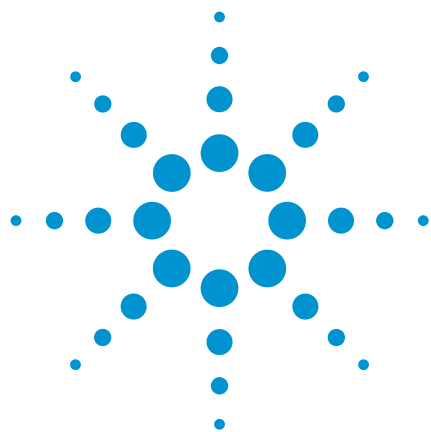




Organochlorine pesticides

Analysis of Arochlor 1016, 1221, 1232 and 1242 to EPA 8080

Application Note

Environmental

Authors

Agilent Technologies, Inc.

Introduction

Gas chromatography with an Agilent CP-Sil 8 CB for Pesticides column separates components in Arochlor 1016, 1221, 1232, and 1242 according to EPA 8080 in 50 minutes.



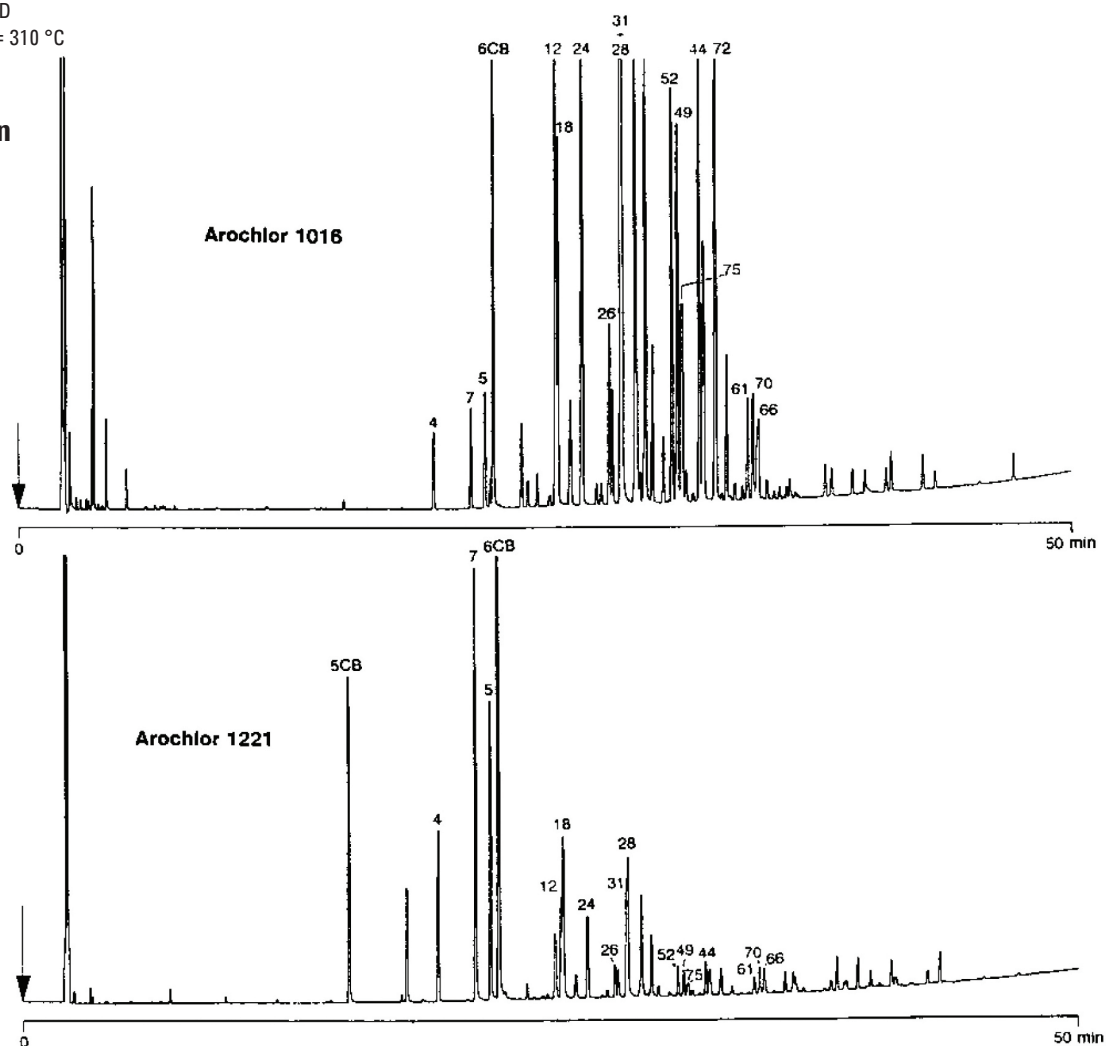
Agilent Technologies

Conditions

Technique : GC-capillary
Column : Agilent CP-Sil 8 CB for Pesticides, 0.32 mm x 50 m
fused silica WCOT CP-Sil 8 CB (df = 0.12 µm)
(Part no. CP7481)
Temperature : 100 °C (2 min) → 275 °C, 3°C/min
Carrier Gas : H₂, 140 kPa (1.4 bar, 20 psi)
Injector : Split, 100 mL/min;
T = 275 °C
Detector : ECD
T = 310 °C

Peak identification

acc. Ballschmider numbers



www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Printed in the USA

31 October, 2011

First published prior to 11 May, 2010

A00797



Agilent Technologies